Benserazide . HCI

Peripheral decarboxylase inhibitor

Peripheral decarboxylase inhibitor. Also inhibits tryptophan hydroxylase and tyrosine hydroxylase. Used with L-DOPA as a Parkinson's disease therapeutic.

Citations: 6

View Online »

Ordering Information

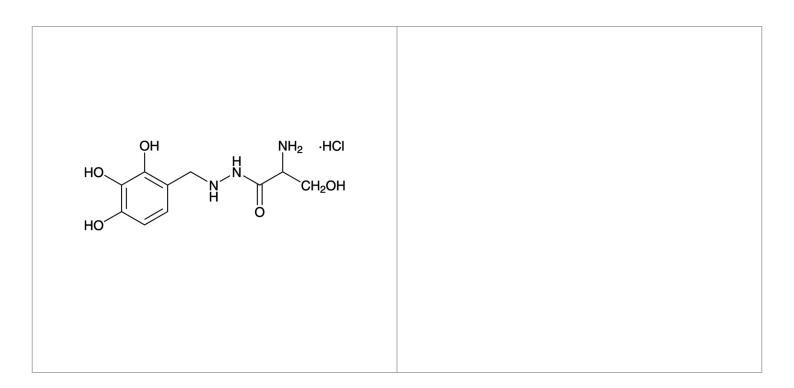
Order Online »

BML-EI165-1000

1g

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Product is subject

to oxidation. Make solutions fresh daily. Store, as supplied, at room temperature for up

to 1 year.

Handling Protect from light.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Tan solid.

CAS 14919-77-8

Couple Target Peripheral decarboxylase, Tryptophan hydroxylase,

Tyrosine hydroxylase

Couple Type Inhibitor

Formula $C_{10}H_{15}N_3O_5$. HCI

MW 293.7

Purity ≥98%

Soluble in water. Almost insoluble in 100% ethanol,

insoluble in acetone.

Last modified: May 29, 2024

